

Product datasheet

Human Factor XI ELISA Kit ab108834

1 图像

概述

产品名称 人Factor XI ELISA试剂盒

检测方法 Colorimetric

精确度

批次内				
样品	n	Mean	SD	CV%
Overall				4.5%

批次间				
样品	n	Mean	SD	CV%
Overall				9.8%

样品类型 Cell culture supernatant, Serum, Plasma

检测类型 Sandwich (quantitative)

灵敏度 = 1.3 ng/ml

范围 6.25 ng/ml - 25 ng/ml

复检 96 %

检测时间 4h 00m

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

产品概述 Abcam's Factor XI Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Factor XI levels in plasma, serum, and cell culture supernatants.

A Factor XI specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a Factor XI specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Complex is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize Streptavidin-Peroxidase enzymatic reaction. TMB is catalyzed by Streptavidin-Peroxidase to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the amount of Factor XI captured in plate.

经测试应用

适用于: Sandwich ELISA

平台

Microplate

性能

存放说明

Store at -20°C. Please refer to protocols.

组件	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent M Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
50X Biotinylated Human Factor XI Antibody	1 x 120µl
Chromogen Substrate	1 x 8ml
Factor XI Microplate (12 x 8 well strips)	1 unit
Factor XI Standard	2 vials
Sealing Tapes	3 units
Stop Solution	1 x 12ml

功能

Factor XI triggers the middle phase of the intrinsic pathway of blood coagulation by activating factor IX.

组织特异性

Isoform 2 is produced by platelets and megakaryocytes but absent from other blood cells.

疾病相关

Factor XI deficiency (FA11D) [MIM:612416]: A hemorrhagic disease characterized by reduced levels and activity of factor XI resulting in moderate bleeding symptoms, usually occurring after trauma or surgery. Patients usually do not present spontaneous bleeding but women can present with menorrhagia. Hemorrhages are usually moderate. Note=The disease is caused by mutations affecting the gene represented in this entry.

序列相似性

Belongs to the peptidase S1 family. Plasma kallikrein subfamily.
Contains 4 apple domains.
Contains 1 peptidase S1 domain.

翻译后修饰

Activated by factor XIIa (or XII), which cleaves each polypeptide after Arg-387 into the light chain, which contains the active site, and the heavy chain, which associates with high molecular weight (HMW) kininogen.

细胞定位

Secreted.

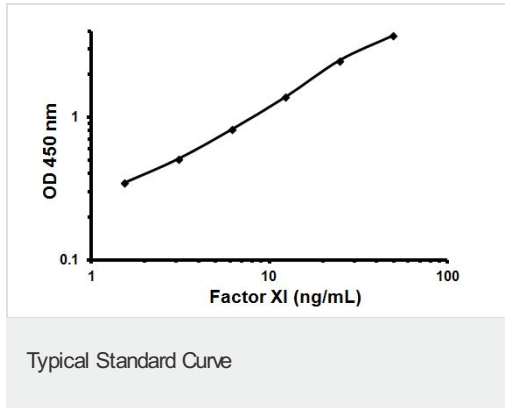
应用

Our [Abpromise guarantee](#) covers the use of **ab108834** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

应用	Ab评论	说明
Sandwich ELISA		Use at an assay dependent dilution.

图片



Representative Standard Curve using ab108834.

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <http://www.abcam.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors